# SVISCISVS

# Product Datasheet

# Cubis® MCM10K3

## Manual Mass Comparator

### User Benefits

- Complete mass standard laboratory in a single unit
- Climate sensors for recording all data relevant for determining measurement uncertainty
- Integrated workflow control for efficient and error-free mass comparison
- Fast measurement cycles according to the ABA, ABBA or AB<sub>1</sub>...B<sub>n</sub>A method

## Highlighted Performance Features

- Cubis® MSA color touch screen for fast and simple configuration of parameters and workflows
- External sensor-equipped climate module for recording the temperature, humidity and air pressure
- Integrated calibration workflows for ABA, ABBA, AB<sub>1</sub>...B<sub>n</sub>A cycles to ensure efficient, error-free mass comparison
- Fully integrated function for determining the measurement uncertainty in accordance with OIML and ASTM recommendations
- Filters for optimal adaptation of the mass comparator to ambient conditions
- Monolithic weighing technology
- For display and evaluation, complete electronics and power supply separated from the weighing system to prevent heat from affecting the results



- This MCM mass comparator features digital eccentric (off-center) load compensation, instead of a moving centering pan, facilitating easier loading of weights
- Additional applications for density determination, statistics and individual identifiers are integrated as standard programs
- Built-in SD card slot for storage and transfer of all data and settings
- Graphical level indicator for interactive user guidance during levelling
- Easy logging of reference weight data
- Continuous weighing range display: any weight between 0 g and the maximum capacity can be displayed
- USB, RS-232C and Ethernet interface ports to integrate the mass comparator into networks or to enable it to communicate with external software via third-party protocols, standardized communication protocols or Web services

# **Technical Specifications**

Metrological Specifications	
Maximum capacity	11 kg
Application range	0 – 11 kg
Readability	1 mg
Repeatability, optimal ')	0.8 mg
Repeatability, standard E <sup>2</sup> )	1 mg
Repeatability, E ½ load ²)	0.8 mg
Repeatability standard, F <sup>3</sup> )	3 mg
Electronic weighing range and tare range	11 kg
Linearity	6 mg
Eccentric load deviation	0.5 mg mm
Stabilization time	3 s
Cycle time, ABBA in s	90 s

Basic Equipment	
Interfaces	RS232C USB LAN
Application programs	Basic weighing, mass unit conversion, individual identifiers, density determination, statistics
Below-comparator weighing port	$\checkmark$
Air temperature sensor	$\checkmark$
Air humidity sensor	$\checkmark$
Air pressure sensor	$\checkmark$
PC connecting cable	USB

Ambient Conditions	
Permissible operating temperature range	10-30 °C
Recommended operating temperature	22 °C
Temperature fluctuations	0.3°C/h 0.5°C/12h
Max. air current	< 0.2 m/s
Humidity range	40-70 %
Humidity fluctuations	5% 4h
Power supply	100-240 V AC/50-60 Hz
Power consumption	< 35 VA

Dimensions	
Weighing pan dimensions ( $W \times D$ )	200 × 200 mm
Sample size (D × H)	240 × 276 × 102 mm
Weigh cell ( $W \times D \times H$ )	239 × 320 × 56 mm
Net weight	9.8 kg
Gross Weight	13.9 kg
Number of packages	1
Package data 1	60 × 40 × 49 cm
Optimal height for setup	800 mm

#### Applications

Applications	
OIML calibration range RS	
OIML calibration range E1	10 kg
OIML calibration range E2	5 kg - 10 kg
OIML calibration range F1	1 kg - 10 kg
OIML calibration range F2	500 g - 10 kg
OIML calibration range M1	200 g - 10 kg
OIML calibration range M2	20 g - 10 kg
OIML calibration range M3	1 g - 10 kg
ASTM E617 calibration range Class 000	
ASTM E617 calibration range Class 00	10 kg
ASTM E617 calibration range Class 0	5 kg - 10 kg
ASTM E617 calibration range Class 1	2 kg - 10 kg
ASTM E617 calibration range Class 2	1 kg - 10 kg
ASTM E617 calibration range Class 3	500 g - 10 kg
ASTM E617 calibration range Class 4	500 g - 10 kg
ASTM E617 calibration range Class 5	100 g - 10 kg
ASTM E617 calibration range Class 6	50 g - 10 kg
ASTM E617 calibration range Class 7	2 g - 10 kg

Optional Accessories	
External calibration weight	10 kg   E2 YCW712-02
Climate module, uncalibrated, for all MCM models	YCM20MC
Calibration of a YCM20MC climate module with DAkkS calibration certificate	YCM20DAkkS
Climate module with DAkkS calibration certificate for all MCM models	YCM20MC-DAkkS
Tower for climate module, for connection to MCM high-capacity models, including cable	YCM20MC-Tower
Optional draft shield	YDS24C
Weighing table	YWT03
Lifting device for 10 kg	YAW51

The standard deviation "s" is the repeatability calculated from 5 ABA cycles under the following conditions:

 Optimal conditions: automatic measurement without operator influence measured in a laboratory under E1 conditions, on a decoupled weighing stone no drafts from above

 Standard conditions, en a decoupled weighing stone in diato non above weighing stone; no drafts from above

3) Standard conditions F: measurement performed mannually in a laboratory under at least F1 conditions, on a non-decoupled weighing stone, air conditioning and minimal drafts from above

#### Germany

Sartorius Lab Instruments GmbH & Co. KG Otto-Brenner-Straße 20 37079 Göttingen Phone +49 551 308 0

### USA

Sartorius Corporation 565 Johnson Avenue Bohemia, NY 11716 Phone +1 631 254 4249 Toll-free +1 800 635 2906

For further information, visit www.sartorius.com